

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP IL-5 Kit
Kit Catalog Numbers: K156QS-Series
Kit Lot Number: K0081001
Expiration Date: 31 Jul 2019

Acceptance Data/ Percent Recovery Specifications:

| Spot | Name | Calibrator | | Cytokine Panel 1 (human) Control 1 | | Cytokine Panel 1 (human) Control 2 | | Cytokine Panel 1 (human) Control 3 | |
|------|----------|------------|-------|------------------------------------|-------|------------------------------------|-------|------------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | NHP IL-5 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|----------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | NHP IL-5 | 6.28 | 80% - 120% | < 20% | 562 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|----------|------------------------------------|------|-------|
| 1 | NHP IL-5 | Calibrator | 844 | pg/mL |
| | | Cytokine Panel 1 (human) Control 1 | 310 | pg/mL |
| | | Cytokine Panel 1 (human) Control 2 | 110 | pg/mL |
| | | Cytokine Panel 1 (human) Control 3 | 34 | pg/mL |

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

Standard Curve Data:

| NHP IL-5 - Spot 1 | | |
|-------------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 844 | 306,000 | 1.7 |
| 211 | 78,600 | 3.0 |
| 52.8 | 19,800 | 2.8 |
| 13.2 | 5,020 | 3.5 |
| 3.30 | 1,310 | 2.7 |
| 0.824 | 374 | 1.5 |
| 0.206 | 141 | 3.1 |
| 0.00 | 64 | 4.9 |

CERTIFICATE OF ANALYSIS

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------------|---------------------|-----------------|
| Human IL-5 Plate | Z0046811 | 2-8°C | 30 Sep 2020 |
| Anti-hu IL-5 Antibody | D0080887 | 2-8°C | 31 Jul 2019 |
| Cytokine Panel 1 (human) Calibrator Blend | A0080121 | 2-8°C | 28 Feb 2022 |
| Diluent 43 | M0430024 | ≤ -10°C | 31 Jan 2020 |
| Diluent 3 | M0030081 | ≤ -10°C | 30 Sep 2020 |
| MSD Read Buffer T | Not Kit Specific | Not Kit Specific | N/A |

| Additional Components Included in V-PLEX Plus | |
|---|------------------|
| Microplate Adhesive Film | Not Kit Specific |
| MSD Wash Buffer (20X), 100 mL | Not Kit Specific |
| Cytokine Panel 1 (human) Control 1 | Not Kit Specific |
| Cytokine Panel 1 (human) Control 2 | Not Kit Specific |
| Cytokine Panel 1 (human) Control 3 | Not Kit Specific |

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 29 MAY 2018 |